

## Intracellular Antibody Delivery with ImmunoFect STAR™ Antibody Transfection Reagent

### Background

ImmunoFect STAR™ is a highly efficient antibody transfection reagent designed for intracellular delivery of antibodies into live cells. ImmunoFect STAR™ enables functional studies of intracellular targets by facilitating direct antibody access to the cytosol. The delivery mechanism is optimized for high efficiency while maintaining cell viability.

### ImmunoFect STAR™ / Antibody Complex Formation

*ImmunoFect STAR™ is supplied in anhydrous ethanol, 10 mg/mL.*

- 1) Dilute with HEPES buffer (20 mM, pH 7.4), (5 µL ImmunoFect STAR™ to 45 µL HEPES buffer), to make a 1 mg/mL ImmunoFect STAR™ solution.
- 2) Prepare the ImmunoFect STAR™ antibody complex by diluting 3 µL of 1 mg/mL ImmunoFect STAR™ with 27 µL HEPES buffer. Combine the 30 µL of diluted ImmunoFect STAR™ solution with 20 µL of antibody solution (2 µL of 1 mg/mL antibody, with 18 µL HEPES buffer) and incubate for 1 minute.
- 3) Add ImmunoFect STAR™ antibody complex solution to 450 µL serum-free DMEM or Leibovitz's L-15 medium and apply to cells.

*It may be necessary to optimize the ratio (ImmunoFect STAR™ concentration of 4-8 µg/mL, and antibody concentration of 2-8 µg/mL) to achieve maximum efficiency for intracellular delivery.*